

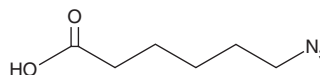
PRODUCT INFORMATION



6-Azidohexanoic Acid

Item No. 16518

CAS Registry No.:	79598-53-1
Formal Name:	6-azido-hexanoic acid
Synonym:	Click Tag™ 6-Azidohexanoic Acid
MF:	C ₆ H ₁₁ N ₃ O ₂
FW:	157.2
Purity:	≥98%
UV/Vis.:	λ _{max} : 214 nm
Supplied as:	A neat oil
Storage:	-20°C
Stability:	≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

6-Azidohexanoic acid is supplied as a neat oil. A stock solution may be made by dissolving the 6-azidohexanoic acid in the solvent of choice, which should be purged with an inert gas. 6-Azidohexanoic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 6-azidohexanoic acid in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 6-azidohexanoic acid can be prepared by directly dissolving the neat oil in aqueous buffers. The solubility of 6-azidohexanoic acid in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

6-Azidohexanoic acid is a six carbon saturated fatty acid with an ω-terminal azide group. This terminal group allows conjugation with compounds containing alkyne groups through a copper(I)-catalyzed cycloaddition reaction, also known as click chemistry. 6-Azidohexanoic acid is commonly used as a linker that is first acylated to one compound to create a derivative that can then be joined to another compound using click chemistry.¹⁻³

References

1. Alleti, R., Rao, V., Xu, L., *et al.* A solanesol-derived scaffold for multimerization of bioactive peptides. *J. Org. Chem.* **75(17)**, 5895-5903 (2010).
2. Damar, P., Wanger-Baumann, C.A., Pillarsetty, N., *et al.* Efficient 18F-labeling of large 37-amino-acid pHLLIP peptide analogues and their biological evaluation. *Bioconjug. Chem.* **23(8)**, 1557-1566 (2012).
3. Zolotarskaya, O.Y., Wagner, A.F., Beckta, J.M., *et al.* Synthesis of water-soluble camptothecin-polyoxetane conjugates via click chemistry. *Mol. Pharm.* **9(11)**, 3403-3408 (2012).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/06/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM